

# Validation data for Anti-HER2-hIgA2

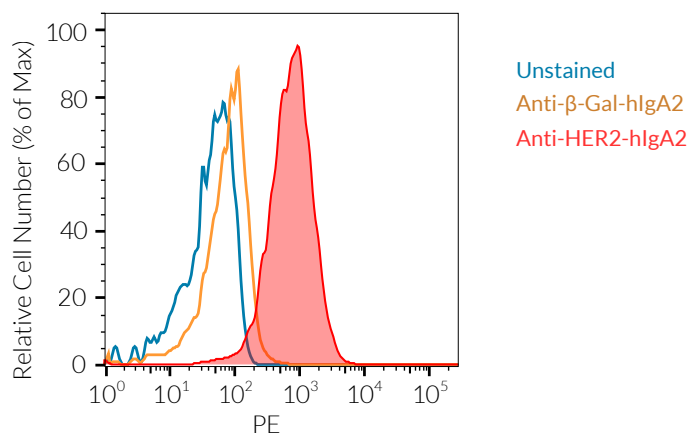
<https://www.invivogen.com/anti-her2-tra-igga2>

For research use only

Version 23K28-AK

Anti-HER2-Tra-hIgA2 is a recombinant monoclonal antibody (mAb) featuring the human epidermal growth factor receptor 2 (HER2)-targeting variable region of the clinical relevant mAb trastuzumab and an hIgA2 constant region. HER2 is found on the surface of epithelial cells and various types of cancer. The detection of HER2 by Anti-HER2-Of-hIgG1 has been verified by flow cytometry using HER2-overexpressing cells (**Figure 1**).

## Binding of Anti-HER2-hIgA2 mAb to HER2



**Figure 1. Binding of Anti-HER2-hIgA2 mAb to HER2.** EL4 cells overexpressing HER2 were incubated for 30 minutes at 4°C with 125 ng of Anti-HER2-hIgA2 or an isotype control. Subsequently, a secondary PE-labeled antibody was added and incubated at 4°C for 30 minutes. The binding affinity was assessed using flow cytometry.

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3-622-34-80

E-mail: [info@invivogen.com](mailto:info@invivogen.com)